

The Advantages of Minisart® Hydrophilic Syringe Filters for HPLC Sample Preparation

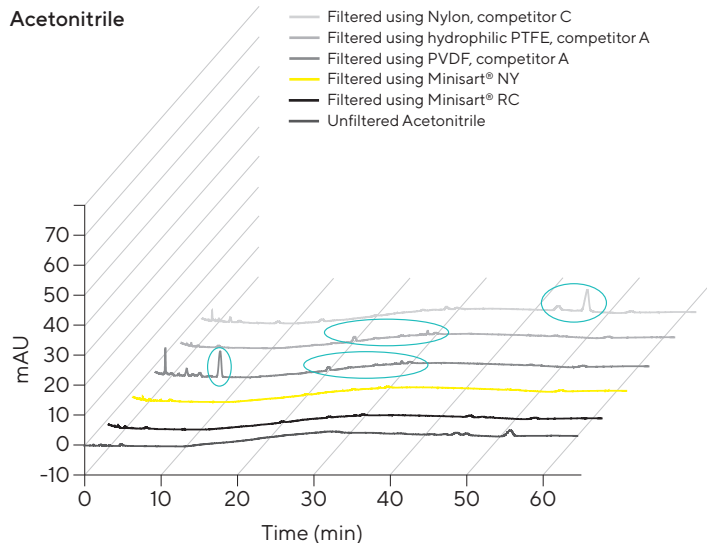
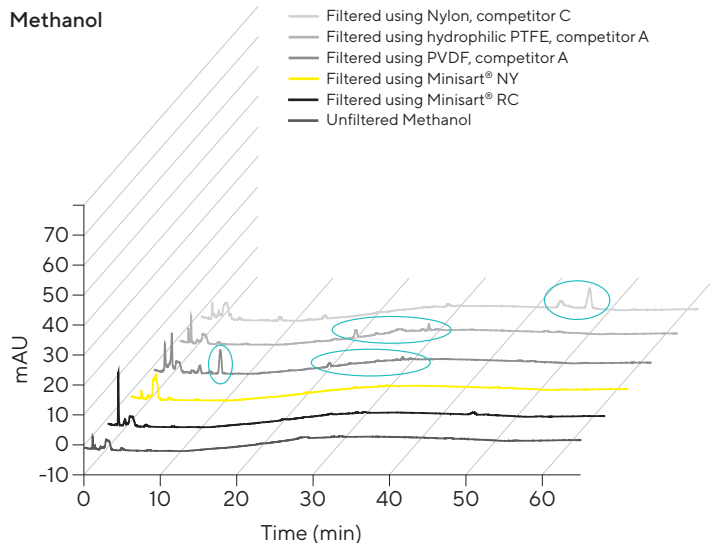
Sartorius Minisart® Versus Other Competitor Hydrophilic Filter Materials

The purpose for using syringe filters for sample preparation prior to analysis is to protect analytical equipment and columns. Sartorius Minisart® filters reliably remove particles from your samples without adding unwanted extractables or leachables (see comparison of different filter membrane materials below). Our re-generated cellulose (RC) membrane used in Minisart® RC and our polyamide (= nylon) membrane incorporated in Minisart® NY are exceptionally pure compared with other common hydrophilic filter membranes and competitor products.

Benefits of Sartorius Minisart®

- Low adsorption of analytes
- Maximum chemical compatibility
- Minimum extractables





HPLC Procedure

Column C18: 250 × 4.6 mm, Flow Rate: 1 mL/min, Wavelength: 210 nm

HPLC Injection Volume: 20 µl, Analysis Time: 65 min, Temperature: 40 °C, Mobile Phases: A) Acetonitrile | B) Water, Gradient: Hold 60% A for 10 min | 60% to 95% A in 20 min | 95% to 100% A in 35 min

Reliable and Ultrapure

Clean & Green

Hydrophilic PVDF and PTFE require an additional coating step with methacrylates to make their membrane surface hydrophilic. Competitor nylon membranes are using raw materials which are especially interfere with LC | MS.

Also, unlike PVDF syringe filters those with RC and nylon membranes are suitable for DMSO and other amides, ketones, esters, and ethers.

Ordering Information

Ø mm	Pore Size	Sterile*	Qty Pk	Order No.
Minisart® RC (Regenerated Cellulose + PP)				
25 mm	0.2 µm	Yes	50	17764-----ACK
25 mm	0.2 µm	No	50	17764-----K
25 mm	0.2 µm	No	200	17764-----S
25 mm	0.2 µm	No	500	17764-----Q
25 mm	0.45 µm	No	50	17765-----K
25 mm	0.45 µm	No	200	17765-----S
25 mm	0.45 µm	No	500	17765-----Q
15 mm	0.2 µm	Yes	50	17761-----ACK
15 mm	0.2 µm	No	50	17761-----K
15 mm	0.2 µm	No	500	17761-----Q
15 mm	0.45 µm	No	50	17762-----K
15 mm	0.45 µm	No	500	17762-----Q
4 mm	0.2 µm	No	50	17821-----K
4 mm	0.2 µm	No	500	17821-----Q
4 mm	0.45 µm	No	50	17822-----K
4 mm	0.45 µm	No	500	17822-----Q

Ø mm	Pore Size	Sterile*	Qty Pk	Order No.
Minisart® NY (Nylon Polyamide + PP)				
25 mm	0.2 µm	Yes	50	17845-----ACK
25 mm	0.2 µm	No	500	17845-----Q
25 mm	0.45 µm	Yes	50	17846-----ACK
25 mm	0.45 µm	No	500	17846-----Q
15 mm	0.2 µm	No	50	1776B-----K
15 mm	0.2 µm	No	500	1776B-----Q
15 mm	0.45 µm	No	50	1776C-----K
15 mm	0.45 µm	No	500	1776C-----Q
25 mm	GF ¹ prefilter 0.2 µm	No	50	1784B-----K
25 mm	GF ¹ prefilter 0.2 µm	No	500	1784B-----Q
25 mm	GF ¹ prefilter 0.45 µm	No	50	1784C-----K
25 mm	GF ¹ prefilter 0.45 µm	No	500	1784C-----Q

* Sterile Minisart® are individually packaged and have been sterilized by ethylene oxide (EO).
 Not presterilized Minisart® can be sterilized by autoclaving at 121 °C for 30 min | or by using EO.
¹ GF prefilter: Ultrapure glass fiber with 0.7 µm particle retention on top of the nylon membrane.

Would You Like to Use Other Membranes?

Please refer to Minisart® SRP (hydrophobic PTFE for highest chemical compatibility).

Are You Looking for CE-certified Minisart®?


Please request CE-certified article numbers for Minisart® NML, Minisart® HY and Minisart® SRP. Sartorius offers a full range of syringe filters dedicated for various filtration applications.

Germany

Sartorius Lab Instruments GmbH & Co. KG
Otto-Brenner-Strasse 20
37079 Goettingen
Phone +49 551 308 0

USA

Sartorius Corporation
565 Johnson Avenue
Bohemia, NY 11716
Phone +1 631 254 4249
Toll-free +1 800 635 2906

 For further information, visit
www.sartorius.com